Caspase-9 monoclonal antibody (10-1-87)

The caspases are a family of cysteine proteases that cleave after certain aspartate residues, and are primarily recognized as mediators of apoptosis. caspases are synthesized as inactive zymogens that can be cleaved to form active enzymes following the induction of apoptosis by stress or death receptors. Initiator caspases (e.g. caspase-8 and -10) are activated by dimerization of the zymogen on a dedicated adaptor protein. These activated initiator caspases in-turn cleave downstream effector or executioner caspases (e.g. caspase-3, -6, and -7) in a cascade-like manner, which cleave key cellular proteins that lead to the morphological changes associated with apoptotic cell death.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

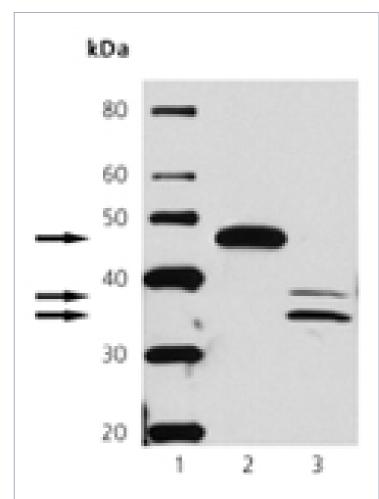
Order Online »

ADI-905-686-100

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of MW marker (1), Jurkat (2), and Jurkat + staurosporine (3) cells probed with Caspase 9 mAb (10-1-87).

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application NotesDetects bands of ~47kDa (full length) and ~37/35kDa

(cleaved) by Western blot.

Clone 10-1-87

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide

and 50% glycerol.

Host Mouse

Immunogen Recombinant full length human Caspase-9.

lsotype lgG2a

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1µg/ml, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Source Purified from mouse ascites.

Species Reactivity Human

UniProt ID P55211

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com